

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT GPR87 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1140

**Lot Number:** C1140-6-090314

**Quantity:** 1 vial ( $2 \times 10^6$ ) frozen cells

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** RH7777

**Transfection:** Expression vector containing full-length human GPR87 cDNA (GenBank accession number NM\_023915.2) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

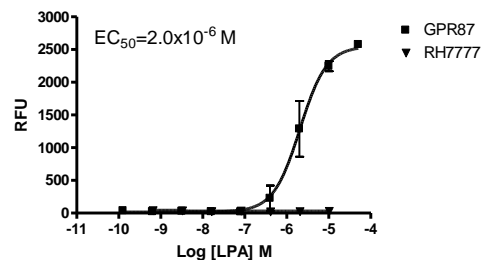
**Propagation Medium:** DMEM, 10% Dialyzed FBS, 3  $\mu$ g/mL puromycin

**Stability:** Stable after minimum two months continuous growth

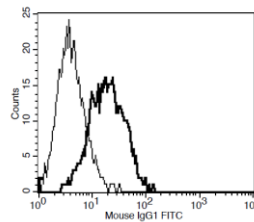
**Background:** GPR87 is a recently discovered LPA receptor. It is highly expressed in neoplastic cells found in the cervical squamous, urinary bladder carcinomas, lung carcinomas, head and neck squamous cell carcinomas, and the placenta. GPR87 plays an important role in normal ovarian development as well as intracellular signal transduction. GPR87 has been linked to lung and ovarian cancer, as well as tumor cell survival. Current research strongly supports the role of GPR87 as a p53 regulator and its potential as a therapeutic target for the treatment of cancer.

**Application:** Functional assays

**Figure 1**



**Figure 2**



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FLIPR. **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

#### References:

Rani, Mukta, et al. "Computational Analysis of the 3-D structure of Human GPR87 Protein: Implications for Structure-Based Drug Design." *Asian Pacific Journal of Cancer Prevention* 14 (2013): 7473-82.

Gugger, Mathias, et al. "GPR87 is an overexpressed G-protein coupled receptor in squamous cell carcinoma of the lung." *Disease Markers* 24 (2008): 41-50.

Zhang, Yanhong, Ariane Scoumanne, and Xinbin Chen. "G Protein-Coupled Receptor 87: a Promising Opportunity for Cancer Drug Discovery." *Mol Cell Pharmacol.* 2 (2010): 111-16.

**FOR RESEARCH USE ONLY.**

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

www.multispaninc.com  
sales@multispaninc.com  
support@multispaninc.com

Ver. October 2005

Phone: +1 (510) 887-0817  
Fax: +1 (510) 887-0863  
26219 Eden Landing Road  
Hayward, CA 94545-3718  
U.S.A.